

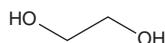
PRODUCT INFORMATION



Ethylene Glycol

Item No. 44724

CAS Registry No.: 107-21-1
Formal Name: 1,2-ethanediol
MF: $C_2H_6O_2$
FW: 62.1
Purity: $\geq 98\%$
Supplied as: A liquid
Storage: Room temperature
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Ethylene glycol is supplied as a liquid. A stock solution may be made by dissolving the ethylene glycol in the solvent of choice, which should be purged with an inert gas. Ethylene glycol is miscible in water.

Description

Ethylene glycol is a cryopreservation agent and synthetic intermediate.^{1,2} It has commonly been used as a major component of antifreeze and coolant.¹ In the lab, ethylene glycol has been used as a cryopreservation agent for various organs and tissues.²⁻⁴ It has also been used as a precursor to polymers and in the synthesis of nanomaterials, among many other uses.¹ This product is also available as part of Cayman's Tissue Cryopreservation Reagent Kit (Item No. 502885).

References

1. Yue, H., Zhao, Y., Ma, X., et al. Ethylene glycol: Properties, synthesis, and applications. *Chem. Soc. Rev.* **41**(11), 4218-4244 (2012).
2. Bojic, S., Murray, A., Bentley, B.L., et al. Winter is coming: The future of cryopreservation. *BMC Biol.* **19**(1), 56 (2021).
3. Watson, R.E., Jr., Wiegand, S.J., Clough, R.W., et al. Use of cryoprotectant to maintain long-term peptide immunoreactivity and tissue morphology. *Peptides* **7**(1), 155-159 (1986).
4. Xu, W., Li, H., and Xu, J. Effective cryopreservation of human brain tissue and neural organoids. *Cell Reports Methods* **4**(5), 100777 (2024).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/06/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM